

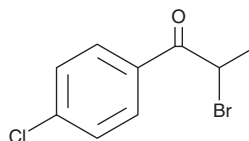
PRODUCT INFORMATION



2-Bromo-4'-chloropropiophenone

Item No. 40861

CAS Registry No.: 877-37-2
Formal Name: 2-bromo-1-(4-chlorophenyl)-1-propanone
Synonyms: 2-Bromo-*para*-chloropropiophenone,
 α -Bromo-4-chloropropiophenone,
 α -Bromo-*para*-chloropropiophenone,
2-Bromo-*p*-chloropropiophenone,
 α -Bromo-*p*-chloropropiophenone
MF: C₉H₈BrClO
FW: 247.5
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

2-Bromo-4'-chloropropiophenone (Item No. 40861) is an analytical reference standard categorized as a precursor in the synthesis of cathinones, including 4-chlorocathinone (Item No. 30379), 4-chloro-N,N-dimethylcathinone (Item No. 24432), 4-chloro-N-isopropylcathinone (Item No. 19392), 4-chloro-N-butylcathinone (Item No. 19972), 4-chloroethcathinone (Item No. 18912), and 4'-chloro- α -pyrrolidinopropiophenone (Item No. 9002179). 2-Bromo-4'-chloropropiophenone is also a precursor in the synthesis of 4-chloromethcathinone (Item No. 16436).¹ This product is intended for research and forensic applications.

Reference

1. Bulska, E., Bachliński, R., Cyrański, M.K., *et al.* Comprehensive protocol for the identification and characterization of new psychoactive substances in the service of law enforcement agencies. *Front. Chem.* **8**, 693 (2020).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/10/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM